



CERTIFICATE OF ANALYSIS

Product Description: Proinflammatory Panel 1 (mouse) Calibrator Blend
Catalog Number: C0048-2
Lot Number: A0080440
Expiration Date: 31 AUG 2030
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	991	pg/mL
IL-1 β	1640	pg/mL
IL-2	2610	pg/mL
IL-4	1710	pg/mL
IL-5	926	pg/mL
IL-6	4950	pg/mL
KC/GRO	2270	pg/mL
IL-10	3440	pg/mL
IL-12p70	31400	pg/mL
TNF- α	602	pg/mL

Reconstitution Volume: 1000 μ L

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Michelle Likiam	Quality		09 OCT 2025